

ReadiLink™ Rapid XFD594 Antibody Labeling Kit

Catalog Number: 1277

Unit Size: 2x50 ug Labelings

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
--------------------	--

Expiration Date	6 months upon receiving
-----------------	-------------------------

Chemical Properties

Appearance	Solid
------------	-------

Molecular Weight	N/A
------------------	-----

Soluble In	DMSO
------------	------

Spectral Properties

Excitation Wavelength	590 nm
-----------------------	--------

Emission Wavelength	618 nm
---------------------	--------

Applications

XFD594 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 594 (Alexa Fluor® is the trademark of Thermo Fisher). Readilink™ protein labeling technology is the most robust and convenient tool for preparing fluorescent antibody conjugates used for fluorescence imaging and flow cytometry applications. ReadiLink™ Rapid XFD594 Antibody Labeling Kit provides the most convenient tool for making XFD594-labeled antibody conjugates. XFD594 dye used in this ReadiLink™ kit is reasonably stable and shows good reactivity and selectivity with protein amino groups. The kit has all the essential components for labeling ~2x50 ug antibody. Each of the two vials of XFD594 dye provided in the kit is optimized for labeling ~50 µg antibody. ReadiLink™ Rapid XFD594 Antibody Labeling Kit requires minimal hands-on time. The prepared XFD594 conjugates (with the kit) are ready to use for fluorescence imaging and flow cytometry applications without further purifications needed. It provides the most convenient method to prepare XFD594-labeled antibodies.